

Human IL-8 ELISA Kit ab46032

★★★★★ 1 Abreviews 62 References 画像数 2

製品の概要

製品名	Human IL-8 ELISA Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Pooled Data	6		CV%
				4%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Pooled Data	6		CV%
				6%
サンプルの種類	Cell culture supernatant, Serum, Plasma			
アッセイタイプ	Sandwich (quantitative)			
検出感度	< 12.3 pg/ml			
検出範囲	31.25 pg/ml - 1000 pg/ml			
添加回収試験	88 %			
全工程の試験時間	2h 15m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Human IL-8 ELISA kit is designed for the quantitative measurement of IL-8 supernatants, buffered solutions, serum and plasma samples. (Genecards: CXCL8)			

A monoclonal antibody specific for IL-8 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-8 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for IL-8 are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of IL-8 present in the samples.

This kit will recognize both endogenous and recombinant Human IL-8.

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

内容	ラベル	2 x 96 tests	1 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7.5ml	1 x 7.5ml
Biotinylated anti-IL-8	Red	2 x 0.4ml	1 x 0.4ml	1 x 0.4ml
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml	1 x 11ml
HRP Diluent	Red	1 x 23ml	1 x 12ml	1 x 12ml
IL-8 Control	Silver	4 vials	2 vials	2 vials
IL-8 Microplate (12 x 8 well strips)		2 units	1 unit	1 unit
IL-8 Standard	Yellow	4 vials	2 vials	2 vials
Plastic Plate Covers		4 units	2 units	2 units
Standard Diluent (Serum)		2 x 7ml	1 x 7ml	1 x 7ml
Stop Reagent	Black	2 x 11ml	1 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl	2 x 5µl

機能

IL-8 is a chemotactic factor that attracts neutrophils, basophils, and T-cells, but not monocytes. It is also involved in neutrophil activation. It is released from several cell types in response to an inflammatory stimulus. IL-8(6-77) has a 5-10-fold higher activity on neutrophil activation, IL-8(5-77) has increased activity on neutrophil activation and IL-8(7-77) has a higher affinity to receptors CXCR1 and CXCR2 as compared to IL-8(1-77), respectively.

配列類似性

Belongs to the intercrine alpha (chemokine CxC) family.

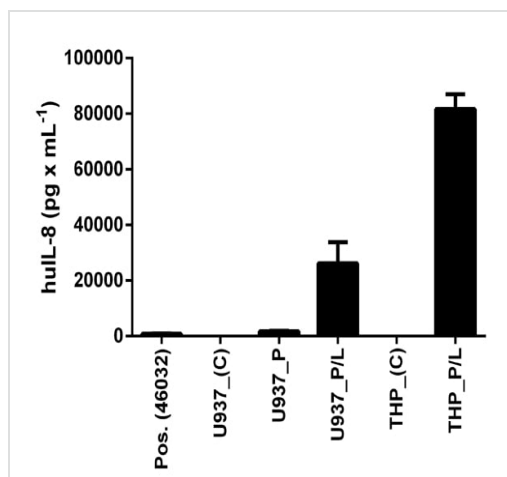
翻訳後修飾

Several N-terminal processed forms are produced by proteolytic cleavage after secretion from at least peripheral blood monocytes, leukocytes and endothelial cells. In general, IL-8(1-77) is referred to as interleukin-8. IL-8(6-77) is the most prominent form.

細胞内局在

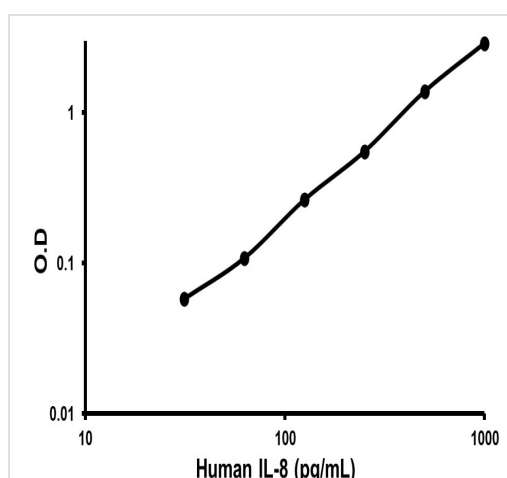
Secreted.

画像



Sandwich ELISA - IL-8 Human ELISA Kit (ab46032)

IL-8 detected in supernatants from control cells (C) or cells stimulated for 24 hours with 50 ng x mL⁻¹ of PMA ([ab120297](#)) (P), and with the addition of 1 ug x mL⁻¹ of LPS (Sigma) (P/L) for the last 6 hours. The positive control is included (Pos.). Results shown after background signal was subtracted (duplicates +/- SD).



Typical Standard Curve

Representative Standard Curve using ab46032

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors